

PRODUCT INFORMATION



BSA- γ -Linolenate Polyunsaturated Fatty Acid Complex (1 mM)

Item No. 39149

Synonym: BSA-GLA
Supplied as: 1 mM γ -linolenate:0.17 mM BSA (6:1 γ -linolenate:BSA) in 150 mM sodium chloride, pH 7.4
Storage: -20°C; avoid freeze/thaw cycles by aliquoting protein
Stability: \geq 1 year
Item Origin: Animal/Bovine

Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Description

BSA- γ -linolenate polyunsaturated fatty acid complex is composed of γ -linolenic acid (Item No. 90220) and bovine serum albumin (BSA) at an approximately 6:1 molar ratio of γ -linolenate:BSA. The BSA/BSA-FAs are acceptable for use with short-term cell culture applications (acute treatment up to 18 hours), however, for long-term applications (25+ hours) the product can/should be filter sterilized using a 0.2 μ m filter and sterile receptacle without affecting the performance. For best results, it is recommended that this product be used in conjunction with BSA control for BSA-fatty acid complexes (1 mM) (Item No. 34932).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 02/11/2025

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897

[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM

WWW.CAYMANCHEM.COM